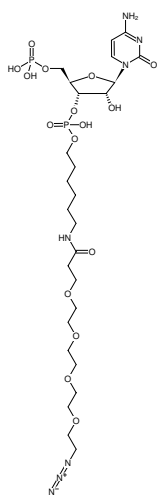




pCp-Azide

Cytidine-5'-phosphate-3'-(15-azido-4,7,10,13-tetraoxa-pentadecanoyl-6-aminohexyl)phosphate, Triethylammonium salt

Cat. No.	Amount
NU-1708	20 µl (1 mM)



Structural formula of pCp-Azide

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₂₆H₄₇N₇O₁₆P₂ (free acid)

Molecular Weight: 775.64 g/mol (free acid)

Exact Mass: 775.26 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Related Products:

pCp-BIO, #NU-1706-BIO

pCp-Desthiobiotin, #NU-1706-Desthiobio

pCp-Cy3, #NU-1706-CY3

pCp-Cy5, #NU-1706-CY5

pCp-Amine, #NU-1706

pCp-Alkyne, #NU-1709

Selected References:

Gerlach *et al.* (2019) Combined quantification of intracellular (phospho-)proteins and transcriptomics from fixed single cells. *Sci. Rep.* **9** (1):1469.